

Nimodipine

Ca²⁺ channel inhibitor

Ca²⁺ channel antagonist. Highly potent at L-type channels. Block some low-voltage activated Ca²⁺ channels. Enhances memory in old or brain-damaged animals. Ameliorates experimental diabetic neuropathy in streptozotocin-induced diabetic rats. Vasodilator.

Citations: 9

[View Online »](#)

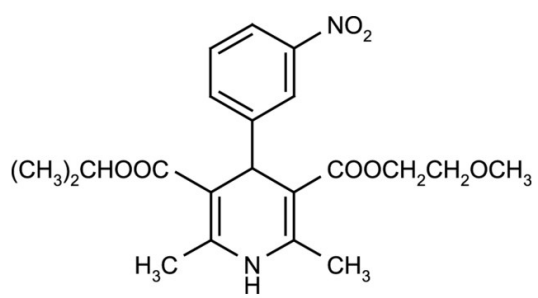
Ordering Information

[Order Online »](#)

ALX-550-277-G001	1g
------------------	----

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability As indicated on product label or CoA when stored as recommended.

Handling Protect from light.

Long Term Storage Ambient

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name 1,4-Dihydro-2,6-dimethyl-4-(3-nitrophenyl)-3,5-pyridinedicarboxylic acid 2-methoxyethyl-1-methylethyl ester

Appearance Yellowish crystalline powder.

CAS 66085-59-4

Couple Target Calcium channel

Couple Type Inhibitor

Formula $C_{21}H_{26}N_2O_7$

MI 14: 6551

MW 418.4

Purity ≥99% (Assay)

Solubility Soluble in DMSO (25mg/ml) or methanol.

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
[info-
eu@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)